

Product Number:	101987
Lot Number:	905916
Expiration:	28-Feb-2020
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.54x10 ⁹ Beads/mL

Review/Approval

Patrizia Stadler

Assoc. Director, QC

p. Stadler 20 Jan 2020

Name

Title

Signature/Date